

Counting Chambers

Product number: CHT4

Lot Number: 2025081301

Reference Bead Batch Number:	25072401	
Reference Bead Qualified Concentration	1.88×10^6 / mL	
Reference Bead Concentration By CHT4	1.92×10^6 / mL	
% Difference	2.2 %	Pass

Specification: Concentration measured by Cellometer™, $C_{\text{cht4}} = \pm 10\%$ of $C_{\text{Test Bead}}$

Manufacturing date: 13-AUG-2025

Document version: D

The information provided in this document is valid for the specified lot number and date of analysis. This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.